

Take  
command  
of your  
gamma  
counting.

revvity



Wizard<sup>2</sup>® Automated Gamma Counters



## The one to count on for years to come

Whether you're conducting research, testing or monitoring, you know that gamma counting delivers unprecedented sensitivity for results you can trust. Now, with the expanded **Wizard<sup>2</sup>** family of automatic gamma counters, you can have outstanding sensitivity together with the highest-level security, connectivity and ease of use, giving you greater command over your gamma counting process. Whatever your application need, there's a Wizard<sup>2</sup> that delivers superior gamma counting performance.

### A gamma counter for today's needs – and tomorrow's challenges

Today's laboratories are pressured to do more with less – fewer resources, less time and tighter turnaround expectations. Wizard<sup>2</sup> gamma counters offer enhanced features to make sure you're up to the challenge now and in the future.

- **Simplicity.** Easy viewing and editing of instrument performance parameters such as protocol definition, conveyor movement, general settings and instrument diagnostic settings
- **Fully automated data analysis.** MyAssays<sup>®</sup> Desktop Pro with QA software for RIA/IRMA and custom data analysis and reporting lets you easily create and transfer custom reports
- **STAT.** Allows high priority, interrupted sample measurement
- **Networkability.** Wizard<sup>2</sup> instruments include Ethernet connections to easily share and analyze your data
- **21 CFR Part 11 compatibility.** With optional MyAssays<sup>®</sup> Desktop Pro Security software, you can automate your audit trail for electronic record and signature compliance
- **More choices than ever.** From routine counting to more sophisticated research to customer validated clinical solutions in a regulated environment, there's a Wizard<sup>2</sup> that's right for your lab
- **Wizard<sup>2</sup> Result Viewer.** Allows data viewing and export of sample data, Instrument Performance Assessment<sup>™</sup> (IPA), normalization and spectral data



# Superior performance labs everywhere depend on

Laboratory professionals have relied on Wizard<sup>®</sup> and Wizard<sup>2</sup> for years. That's because Wizard<sup>2</sup> instruments have consistently provided superior gamma counting performance. Only Wizard<sup>2</sup> counters provide the following high-performance features.

## Well-type detectors for superior counting efficiency

Well-type NaI (Tl) detectors collect signals from all around the sample rather than from the sides only. This minimizes the variance effect caused by sample position and volume and provides the best counting efficiency possible.



Well-type detector.

## Effective lead shielding for consistent background and minimal crosstalk

In a universal gamma counter, the lead shielding around the detector is another important factor in determining the performance of the instrument. Wizard<sup>2</sup> counters feature a unique sample deck, well type detectors and lead shielding to decrease interferences such as detector and sample deck crosstalk, environmental and background radiation. Additionally, they feature the most lead shielding available on gamma counters today. As a result, you can rely on Wizard<sup>2</sup> counters to surpass all other gamma counters in background and crosstalk reduction.

Wizard<sup>2</sup> Lead Shielding Measurements:

Instruments	Detector Lead Thickness (mm/in)	Lead Shielding Thickness (mm/in)	Weight (kg)
2470 Wizard <sup>2</sup>	30/1.2	12/0.5	57
2480 Wizard <sup>2</sup>	75/3.0	50/2.0	210



# Efficiency that delivers

## Robotic sample changer further minimizes crosstalk

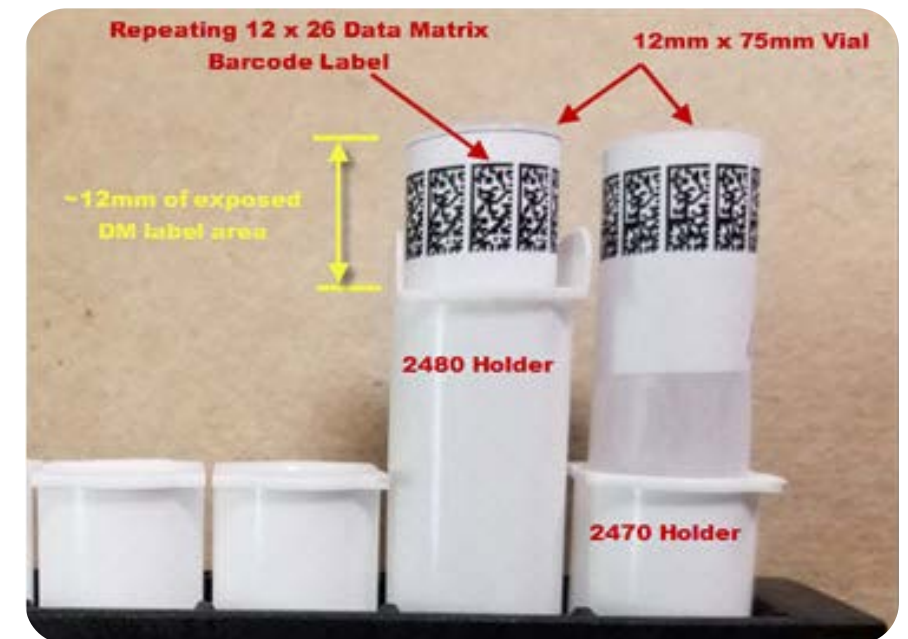
Only Wizard<sup>2</sup> counters feature a unique sample changer system that enables the lifting of samples to the well-type detectors from a conveyor. The conveyor, with the remaining samples, always remains hidden behind the lead shielding.



Robotic sample changer.

## Positive ID sample barcode

The available Sample Vial Barcode Option provides an easy way for the Wizard<sup>2</sup> instrument to positively identify each sample vial using a barcode label attached to the top of each vial.



Example sample ID barcodes.

# MyAssays Desktop

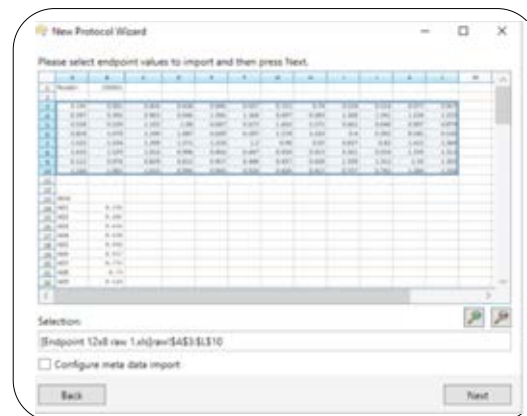


First-class data analysis tools for bioassays

## Easy to start

MyAssays<sup>®</sup> Desktop will kick-start your analysis with preconfigured protocols, quick start wizards and accompanying online content to get you up and running faster than ever.

- Ready-to-use and easy-to-edit preinstalled example protocols
- Wizards guide you through building your own tailor-made protocols
- Portable protocol format means you can easily share and re-use methods with colleagues, from tech-support and online protocols ([www.myassays.com](http://www.myassays.com))
- Intuitive layout editor designed for rack-based assays (easily arrange, and change your sample mapping)
- Integrated seamlessly with Wizard<sup>2</sup> counting

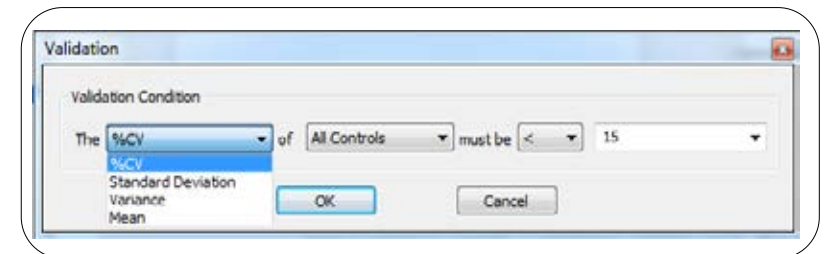


| Specifying intra-assay validation rules.

## Ready for performance

All the data analysis tools you would expect and a whole lot more:

- Transform your data just as your assay dictates: background/blank corrected, % binding, B/B0%, B/T%, factor based, etc.
- Quantitative tools: weighted and unweighted 4-parameter-logistic, 5-parameter-logistic, linear regression, cubic spline, cubic regression, point-to-point and more
- Qualitative tools: automatically label your sample results with cut-off points (and optional grey zones)
- Easy layout-linked replicate analysis (%CV, standard deviation, variance, median, average, etc.)
- Validation: easy to configure intra-assay and inter-assay validation rules. Flag, highlight or add warning messages to your reports
- Detect and mark outliers manually, automatically or directly from the curve to exclude from calculations



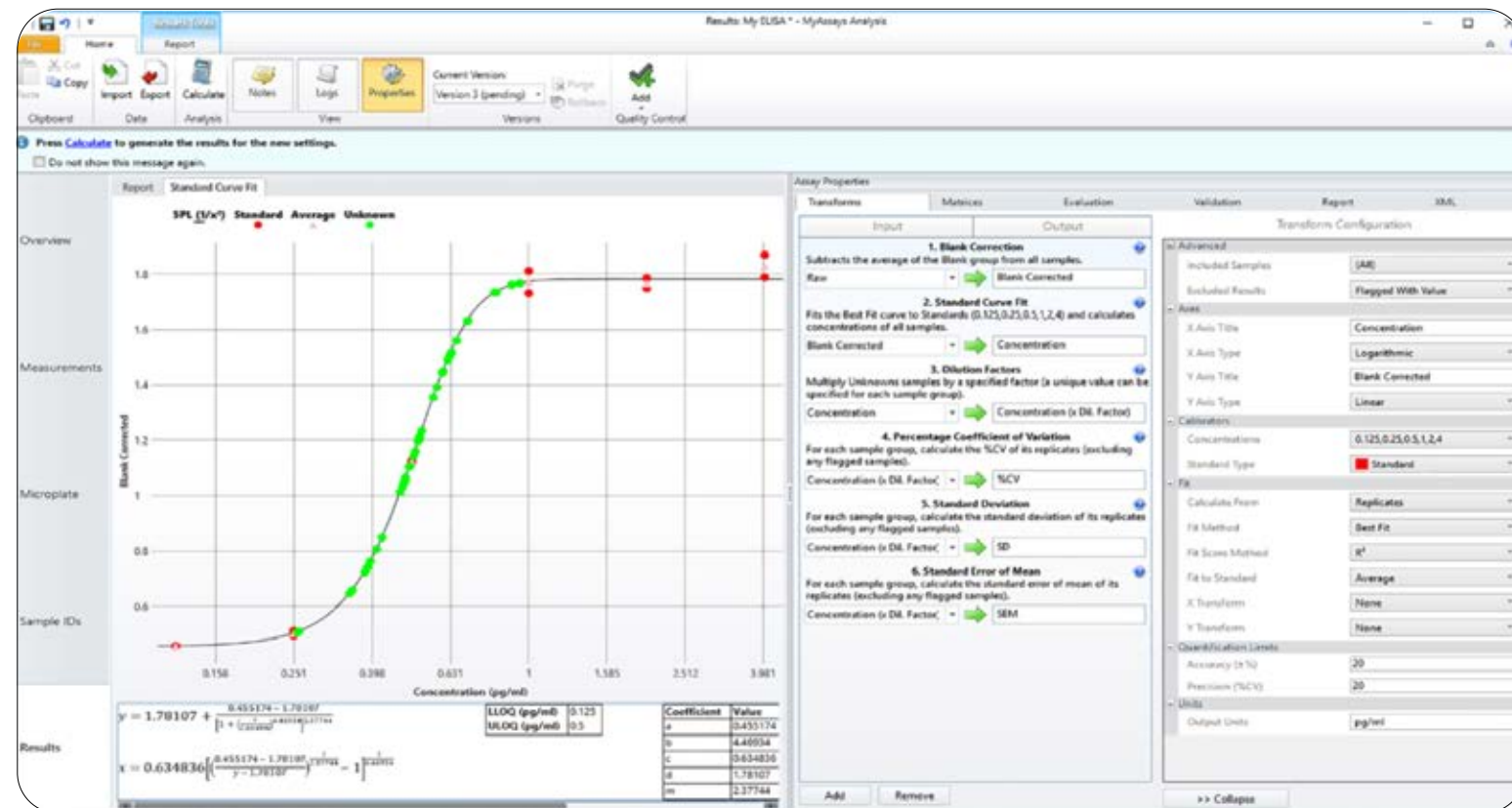
| Specifying validation condition.



### Power and flexibility

MyAssays<sup>®</sup> Desktop data analysis framework neatly handles more sophisticated applications:

- Create your own methods on top of the included extendable library of data-analysis components designed for assays
- Use expressions to write your own formula (refer directly to sample measurements, groups, calculated parameters and reduced data)
- Flexible reporting options - export to Excel/Word/HTML/plain text or auto-print
- Worklists functions - automatically import sample IDs, concentrations, other data analysis parameters, and assay data fields into analyses and reports
- Advanced import - utilize preconfigured and customizable import scripts to get your data into whatever text/CSV file format you use
- Advanced export options designed to produce exacting output compatible with your LIM system



Ready-to-run example protocols in a familiar user-interface.



### Quality control

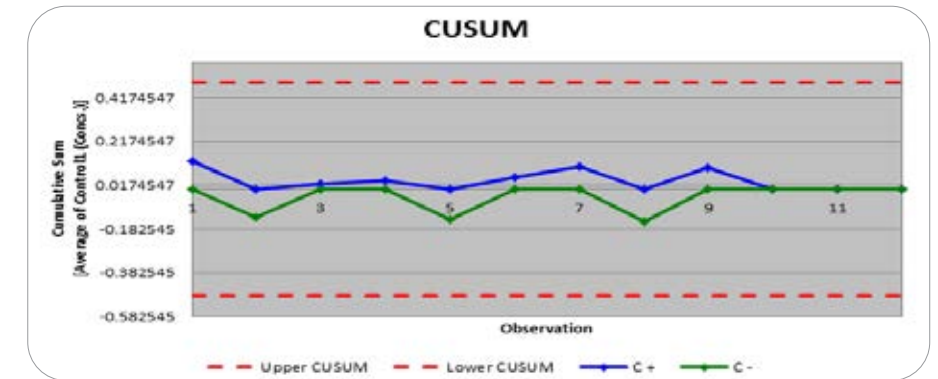
Process inter-assay observations for charting, monitoring and validation of real-time and/or historical data

- Flexibly define multiple observations to track (such as control measurements, analysed data, ED20, ED50, ED80, etc.)
- Generate CUSUM, Levey-Jennings charts as PDF or Excel
- Add automatic warning messages to assay reports to automatically report inter-assay QC violations
- Observe on-demand or as part of a MyAssays<sup>®</sup> Desktop protocol

### 21 CFR Part 11 compatibility

Optional Enhanced Security (ES) module designed specifically for integration with Wizard<sup>2</sup> system provides features required for 21 CFR Part 11:

- Data files protected at the operating-system level for maximum security against unauthorized access
- Shared authentication system with Wizard<sup>2</sup> Enhanced Security system to simplify user administration
- User management – assign each user to a customizable group of authorizations
- Audit trail – secured exportable and printable logs containing full details of all log-in attempts, user operations and full calculation history all date/time stamped and user linked



CUSUM chart observing inter-assay control concentrations.

### MyAssays<sup>®</sup> Desktop Online Support

- Renowned technical support experts.
- Extensive online training videos provided at [www.myassays.com](http://www.myassays.com)
- Further assay protocols available online or through technical support



# Which Wizard<sup>2</sup> is right for you?

Wizard<sup>2</sup> gamma counters offer you choice in sample capacity, shielding thickness and configurations, including optional security features. With so many options, there's a Wizard<sup>2</sup> that's right for you and your applications. No matter which model you choose, you can count on superior gamma counting performance year after year.

## For research applications

### 2470 Wizard<sup>2</sup> with optional MyAssays® Desktop data analysis

- **Ideal for RIA and IRMA studies.** All RIA tube-based studies can be performed with Wizard<sup>2</sup> instruments. Optional MyAssays® Desktop software can be added for additional data analysis
- **Perfect for chromium release studies.** Minimal crosstalk from samples on the conveyor means that the Wizard<sup>2</sup> is optimal for working with higher energy isotopes such as <sup>51</sup>Cr
- **Isotope library.** Contains information for 45 radionuclides
- **IPA.** Wizard<sup>2</sup> monitors 9 detector parameters and automatically provides documentation
- **LAN connectivity and USB.** Makes networking and data transfer flexible
- **Choice of models.** Available with 1, 2, 5 or 10 detectors with 550-sample capacity; 5 and 10 detectors with 1,000-sample capacity
- **Compact footprint.** The 550-sample Wizard<sup>2</sup> is the smallest automatic 10-detector gamma counter available. Its 65 x 77 cm (25.6 x 30.3 in) footprint makes the most of your lab space



Wizard<sup>2</sup> is available in several models featuring 550- and 1,000-sample configurations and with 1, 2, 5 or 10 detectors.





# Which Wizard<sup>2</sup> is right for you?

For research, analytical or customer clinical applications

## 2480 Wizard<sup>2</sup> with optional MyAssays® Desktop ES (21 CFR for PET compatibility)

- **Superior performance for PET studies.** Designed for minimal crosstalk, the 2480 Wizard<sup>2</sup> delivers the best possible counting performance for PET studies
- **Ideal for high-energy or low-activity samples.** With its high-efficiency detector, ability to count sample volumes up to 20 mL and effective lead shielding, the 2480 Wizard<sup>2</sup> is particularly suitable for high-energy gamma emitters, low-activity samples, environmental samples or Schilling tests
- **High-efficiency, 3-inch well-type detector.** Delivers improved counting efficiency, extended energy range and better energy resolution
- **Isotope library.** Contains information for 51 radionuclides
- **One model for all special application needs.** Accommodates both 1,000- and 270-sample capacity of 13 mm (0.5 in) and 28 mm (1.1 in) vial diameters, respectively
- **IPA.** Wizard<sup>2</sup> monitors 9 detector parameters and automatically provides documentation
- **LAN connectivity and USB.** Makes networking and data transfer easy

What's more, both models deliver Enhanced Security compatibility and LIM system integration, with optional MyAssays® Desktop ES (21 CFR compatibility) and Wizard<sup>2</sup> Enhanced Security Software to streamline 21 CFR part 11 compatibility and your ability to upload work lists and download results to a LIM system.



Wizard<sup>2</sup> is available in several models featuring 550- and 1,000-sample configurations and with 1, 2, 5 or 10 detectors.



## Superior performance, yet easy to use

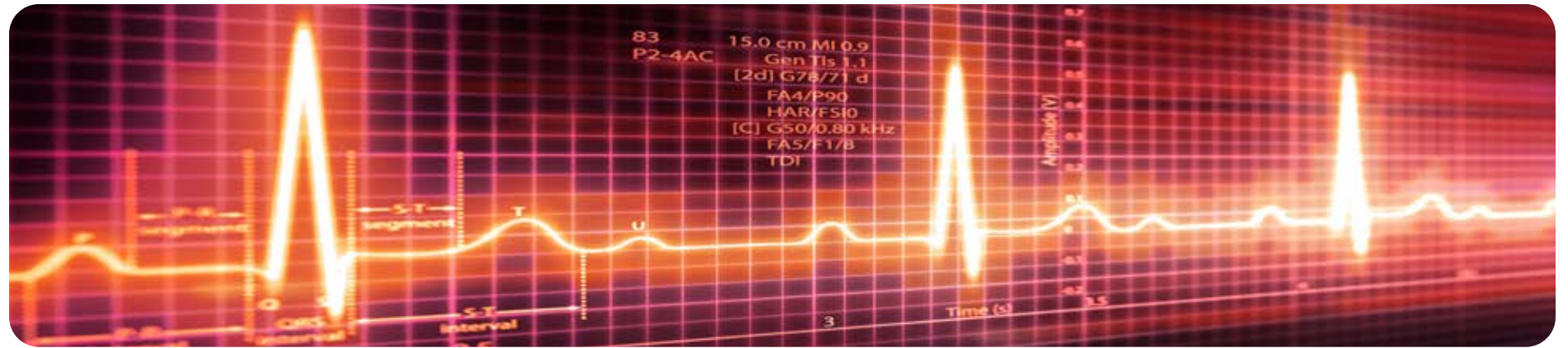
No matter which model you choose, Wizard<sup>2</sup> is designed for ease of use. Once protocols are entered, simply load your samples and press **START**. It's that simple to take command of your gamma counting.

### Revvity IPA: Designed to detect any errors before you do

Exclusive to Wizard<sup>2</sup>, IPA automatically supports laboratory QC systems by performing routine tasks for continued reliability.

- Instrument Performance Assessment (IPA) inspects the isotope spectrum and immediately warns of any condition that generates an unexpected result
- IPA automatically monitors parameters such as window counts and absolute efficiency on each detector and provides documentation
- IPA offers at-a-glance performance history views on the simple touch screen, which further improves your lab's productivity





**Wizard<sup>2</sup>: Choose the model and options right for your labs**

	2470 Wizard <sup>2</sup>		2480 Wizard <sup>2</sup>
Sample capacity available	550	1,000	270/1,000*
Number of detectors available	1, 2, 5 or 10	5 or 10	single 3-inch detector
Wizard <sup>2</sup> Data Viewer	optional	optional	standard
Wizard <sup>2</sup> Enhanced Security	optional	optional	optional
MyAssays <sup>®</sup> Desktop	optional	optional	optional
MyAssays <sup>®</sup> Desktop ES (21 CFR compatibility)	optional	optional	optional
Positive Sample Barcode Reader	optional	optional	optional

\* Model can run 270 20 mL vials or 1,000 12 x 75 mm tubes

**Count on Wizard<sup>2</sup>**

From setup to training, low to high capacity, simple counting studies to sophisticated analysis, look to Wizard<sup>2</sup> as the long-term solution that delivers superior counting performance and excellent security. It's the one your lab can count on for years to come.



## Results. Smarter. Faster.

With a growing emphasis on translational insight, it is more important than ever to be able to examine the molecular mechanisms of disease and translate your *in vitro* models into *in vivo* results. Revvity offers leading solutions and renowned expertise in assays, imaging and Revvity Signals that will help you bring it all together. Whether working in a well, cells or small animals, now you can focus on your science, gain insight sooner and succeed faster.



[www.revivity.com](http://www.revivity.com)

revivity

Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
[www.revivity.com](http://www.revivity.com)

For a complete listing of our global offices, visit [www.revivity.com](http://www.revivity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.

1117301 (383321)